



RFE MEMO


Page 1 of 1

Product Description(s): MSD GOLD™ Read Buffer

Catalog Number: R92TG-Series

Lot Number: Y0600006

Based on additional testing, the original date of expiration for this buffer has been extended by 12 months. The expiration date of Y0600006 has been extended to 30 NOV 2022.

Function	Name	Signature	Date
Quality	Karen Hamilla		20 OCT 2021

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ Read Buffer

Catalog Number/Size: R92TG-1

Lot Number: Y0600006

Expiration Date: 30 NOV 2021

Storage Temperature: Room Temperature

Statement: MSD GOLD Read Buffer is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of MSD GOLD Read Buffer are provided in the product insert available at www.mesoscale.com.

MSD GOLD Read Buffer was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010605 and MSD GOLD 96-well Streptavidin SECTOR® Plate lot # Z0021294.

Concentration of BTI	Mean ECL spec and tolerance range	Intraplate CV or SD spec	Pass/Fail?
2.0 nM (0.10 pmol)	25,531 counts \pm 15%	10 % CV	PASS
0.05 nM (0.0025 pmol)	525 counts \pm 15%	10 % CV	PASS
0 nM (0 pmol)	58 \pm 10 counts	15 count SD	PASS

Statement: This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Julita Arriaga	Quality Control	<i>J. Arriaga</i>	13 DEC 2018
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	13 DEC 2018